

Custom CRISPR Application Form

Instructions:

The information provided will be used to recommend designs for your project. Please answer all questions in the spaces provided to help expedite your request. After completing the form, please email to CRISPR@sial.com.

For an example of a completed CRISPR Application, please [click here](#).

Date

1. Target Sequence	
1a. What genome are you interested in editing? (Please select one) <input type="checkbox"/> Human <input type="checkbox"/> Mouse <input type="checkbox"/> Rat <input type="checkbox"/> Other (genus species):	
1b. Name of target gene and/or Gene ID:	1c. NCBI Ref Seq Number (e.g. NM_000111): * Need help finding Ref Seq? Click here .
1d. Please paste target genomic DNA sequence below:	
1e. Please specify additional transcripts to be targeted below:	
2. Project Goals	
2a. Please indicate the type of project:	
<input type="checkbox"/> Gene Knockout <input type="checkbox"/> Gene Modification (Please see options below. Use Section 7 if necessary.)	
Conditional Knockout <i>*Please paste sequence to be deleted.</i> <i>*See Section 5 for LoxP information.</i>	
SNP Modification <i>*Please specify 25 nt of genomic sequence upstream and downstream with modification(s) highlighted. (e.g. A>T)</i> <i>Example:</i> <i>Before: ACACCCTTCACGTCAGCAGTCGGCACTG</i> <i>AGTCCCTGGCACCAGGCTCTGG</i> <i>After: ACACCCTTCACGTCAGCAGTCGGCICTG</i> <i>AGTCCCTGGCACCAGGCTCTGG</i>	Before:
	After:

Insertion <i>*Please paste insertion sequence and 25 nt upstream and downstream of insertion site.</i> <i>Example:</i> ACACCC TTCACGTCAGCAGTCGGCA_C TGAGTCCCT GGCACCAGGCTCTGG	Insertion Sequence:		
	Insertion Site:		
Deletion <i>*Please paste sequence to be deleted.</i>	Deletion Sequence:		
2b. What is the model system for your project? Cell Line and/or Whole Animal			
3. CRISPR Format Please indicate the format(s) you would like for your CRISPRs:			
3a. Plasmid			
All-in-One Vector gRNA + Cas9 WT	GFP	or	RFP
Dual Vector Cas9 gRNA	GFP	or	RFP or No Fluorophore
Paired Nickase Cas9D10A Paired gRNA	GFP	or	RFP or No Fluorophore
3b. RNA and Protein			
gRNA	Synthetic gRNA	or	<i>In vitro</i> transcribed gRNA
Cas9 mRNA			
Cas9D10A Nickase mRNA			
Cas9 Protein			
3c. Lentivirus			
Virus Particles			
Plasmid			
Vector:	Cas9 Only Options Cas9 Neo <i>(EF1a-Cas9-2A-Neomycin)</i> Cas9 Blast <i>(EF1a-Cas9-2A-Blasticidin)</i>	gRNA Only Options Gecko Puro <i>(U6-gRNA/human-teF1a-puro)</i> Gecko eGFP <i>(U6-gRNA/human-teF1a-tGFP)</i>	All-in-One Options All-in-One <i>(U6-gRNA/truncated human-eF1a-puro-2A-Cas9-tGFP)</i> Sanger <i>(U6-gRNA/hPGK-puro-2A-tBFP)</i>
3d. CRISPR Activator		Activation Region Sequence:	
dCas9p300 Activator gRNA <i>(10 tiled gRNAs)</i> <i>*Please paste activation region sequence.</i>			

4. Validation		
4a. Would you like to order Sigma-Aldrich CRISPR Validation Service to test gRNA cleavage efficiency?		Yes No
5. Donor Design		
5a. Would you like to order Donor Design Service?		Yes No
5b. Preferred Donor Format:		ssDNA oligo or dsDNA plasmid
5c. Conditional KO LoxP Sequence Information <i>Sequences below will be used unless alternate sequences are provided.</i>		
Lox P Upstream ATAACTTCGTATAATGTATGCTATACGAAGTTAT		
Lox P Downstream ATAACTTCGTATAATGTATGCTATACGAAGTTAT		
6. Contact Information		
6a. Title:	6b. First Name:	6c. Last Name:
6d. Institution or Company Name:		
6e. Shipping Address (building name and street address):		
6f. City/Town, State/Province, and Country (include ZIP code or postal code where appropriate):		
6g. Email Address:	6h. Telephone (include country and area code):	
6i. Name of Sales Representative:		
6j. May we contact regarding future promotions?		Yes No
7. Additional Information		
7a. Please provide any additional project details below:		

Merck KGaA
Frankfurter Strasse 250
64293 Darmstadt, Germany

To place an order or receive technical assistance

Order/Customer Service: SigmaAldrich.com/order
 Technical Service: SigmaAldrich.com/techservice
 Safety-related Information: SigmaAldrich.com/safetycenter

Web Order: SigmaAldrich.com/CRISPR

